

ProLong[®] Diamond Antifade Mountant**ProLong[®] Diamond Antifade Mountant with DAPI**

Pub. no. MAN0010261

Revision B.0

Detailed protocol is available online at www.lifetechnologies.com/manuals.

Product	Amount	Cat. no.	Storage*
ProLong [®] Diamond Antifade Mountant	5 × 2 mL 10 mL	P36961 P36970	• 2–8°C • Protect from light
ProLong [®] Diamond Antifade Mountant with DAPI	5 × 2 mL 10 mL	P36962 P36971	

Refractive index of cured mountant: 1.47

* Product may also be stored at $\leq -20^{\circ}\text{C}$. When stored as directed, the product is stable for at least 6 months.



1. Add 1 drop of ProLong[®] Diamond Antifade Mountant onto glass slide.



2. Remove excess liquid from coverslip.



3. Place coverslip sample-side down onto ProLong[®] Diamond Antifade Mountant on the glass slide.



4. Cure for 24 hours at room temperature in the dark.



5. Clean coverslip and image cells.

Limited Product Warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

DISCLAIMER - LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

©2014 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

For support visit lifetechnologies.com/support
lifetechnologies.com